
KLRG1 Polyclonal Antibody

(Catalog #A55655)

Background

Plays an inhibitory role on natural killer (NK) cells and T-cell functions upon binding to their non-MHC ligands. May mediate missing self-recognition by binding to a highly conserved site on classical cadherins, enabling it to monitor expression of E-cadherin/CDH1, N-cadherin/CDH2 and R-cadherin/CDH4 on target cells.

Description

KLRG1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Preservative: 0.02% sodium azide Constituents: 50% Glycerol, PBS, pH 7.3

Specificity

Human

Isotype

IgG

Uniprot ID

Q96E93

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Killer cell lectin-like receptor subfamily G member 1 protein (60-195AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

C-type lectin domain family 15 member A, ITIM-containing receptor MAFA-L, MAFA-like receptor, Mast cell function-associated antigen, KLRG1, CLEC15A, MAFA, MAFAL

Application

ELISA, IHC, IF; Recommended dilution: IHC 1:20 - 1:200, IF 1:10 – 1:100
